



SPECIFICATIONS

CREOPTIX WAVEdelta

GENERAL	
Noise (RMS)	<0.01 pg/mm ² @ 1 Hz
Drift	<0.3 pg/mm ² /min
Readout Frequency	1 Hz, 10 Hz or 40 Hz
Association Const. Range	$k_a = 10^3 - 5 \times 10^7 \text{ M}^{-1} \text{ s}^{-1}$ (small molecules) $k_a = 10^3 - 3 \times 10^9 \text{ M}^{-1} \text{ s}^{-1}$ (large molecules)
Dissociation Const. Range	$k_d = 10^{-5} - 10 \text{ s}^{-1}$
Analysis Temperature Range	4°C - 45°C (max 20°C below ambient)
Molecular Weight Limit	No lower limit
waveRAPID® Functionality	Yes
FLUIDICS	
Flow Channels / Path	4, parallel
Channel Referencing	Any combination of the 4 channels
Flow Cells	Sealed, disposable, integrated into disposable WAVEchip
Flow Rate	1 - 400 µl/min
Crude Sample Robustness	Yes
SAMPLE HANDLING	
Sample Capacity	2x microtiter plates (96 or 384 well, standard or deep well) or vial racks (48 positions of 1.5ml)
Buffer	Automatic switching between 4 buffers
Degasser	Built-in
Injection Volume	< 450 µl, 100 µl typical
Sample Volume Required	Injection volume plus 15-50 µl (application dependent)
Sample Storage Temperature	Ambient or 4°C - 20°C regulated
Sample Recovery	Yes
Automation	120h of unattended operation
DATA TREATMENT	
Information Provided	Kinetic and affinity data (k_a , k_d , K_D)
Graphs	Real-time curves, multiple curve overlays, fit, report point plots
Data Extraction	Curves, k_a , k_d , K_D tables, graphs, reports
Data Analysis	Fully automated data evaluation
Kinetic Models	Predefined models including 1:1 interaction, mass transport, heterogenous ligand, conformational change and bivalent
Direct Kinetics	Yes